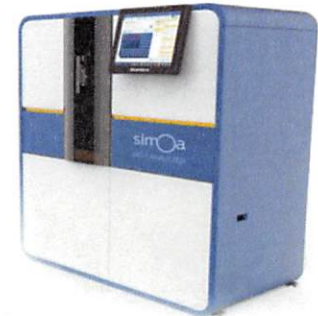


Product Number:	102504
Lot Number:	501195
Expiration:	28-Sep-2018
Platform(s):	HD-1



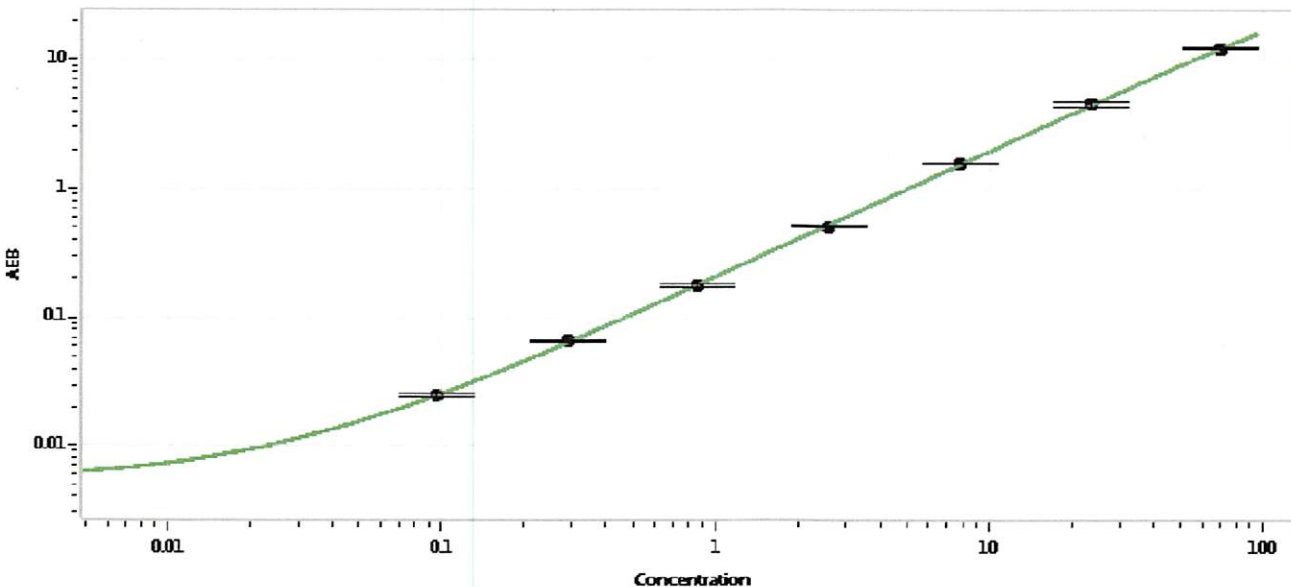
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102497	MF-040218-05
Detector Stock	102498	MF-040218-02
Calibrator Concentrate	102501	MF-022817-06
Calibrator Diluent	102499	MF-032818-09
Sample Diluent	102500	MF-040318-04

Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Concentrate		101957
RGP Reagent		103159

Calibrator Concentrate Target Value	1.4 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.096
C	0.288
D	0.864
E	2.59
F	7.77
G	23.3
H	70.0

Review/Approval: Sivasankari Murugan Name, QC Supervisor Title, [Signature] Signature/Date, 12 APR 2018 Date